## Guanine nucleotide-binding protein Alpha-q / GNAQ / G-Alpha-q Monoclonal Mouse Antibody (GNAQ/2434)



## **Product Description**

GNAQis 359 amino acids long and is identical in all but 1 amino acid residue to the mouse protein. Analysis of human genomic DNA revealed an intronless sequence with strong homology to human GNAQ cDNA. In comparison to GNAQ cDNA, this genomic DNA sequence included several small deletions and insertions that altered the reading frame, multiple single based changes, and a premature termination codon in the open reading frame, all hallmarks of a processed pseudogene. Probes derived from human GNAQ cDNA sequence mapped both chromosomes 2 and 9 in higher constringency genomic blot analyses of DNA from a panel of human/rodent hybrid cell lines.

Primary antibodies are available purified, or with a selection of fluorescent CF® dyes and other labels. CF® dyes offer exceptional brightness and photostability. See the CF® Dye Brochure for more information. Note: Conjugates of blue fluorescent dyes like CF®405S and CF®405M are not recommended for detecting low abundance targets, because blue dyes have lower fluorescence and can give higher non-specific background than other dye colors.

**Stock status:** Because Biotium offers a large number of antibody and conjugation options, primary antibody conjugates may be made to order. Typical lead times are up to one week for CF® dye and biotin conjugates, and up to 2-3 weeks for fluorescent protein and enzyme conjugates. Please email <a href="mailto:order@biotium.com">order@biotium.com</a> to inquire about stock status and lead times before placing your order.

Catalog number key for antibody number 2434, Anti-G-Alpha-q|GNAQ|Guanine nucleotide-binding protein Alpha-q (GNAQ/2434)

## Product attributes

Call us: 800-304-5357

Product attributes				
Antibody number	#2434			
Antibody reactivity (target)	G-Alpha-q, GNAQ, Guanine nucleotide-binding protein Alpha-q			
Antibody type	Primary			
Host species	Mouse			
Clonality	Monoclonal			
Clone	GNAQ/2434			
Isotype	lgG2b, kappa			
Molecular weight	45 kDa			
Synonyms	CMC1; G alpha Q; G protein alpha Q; GAQ; guanine nucleotide binding protein (G protein); q polypeptide; Guanine nucleotide-binding protein G(q) subunit alpha; SWS			
Human gene symbol	GNAQ			
Entrez gene ID	2776			
SwissProt	P50148			
Unigene	269782			
Immunogen	Recombinant full-length human GNAQ protein			
Antibody target cellular localization	Plasma membrane			
Species reactivity	Human			
Species reactivity  Antibody application notes	Human  For coating for ELISA, order Ab without BSA, Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Optimal dilution and staining procedure for a specific application should be determined by user, Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry			
	For coating for ELISA, order Ab without BSA, Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Optimal dilution and staining procedure for a specific application should be determined by user, Recommended starting concentrations for titration are 1-2 ug/mL for most			
Antibody application notes	For coating for ELISA, order Ab without BSA, Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Optimal dilution and staining procedure for a specific application should be determined by user, Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry  HL-60, MOLT-4 and MCF-7 cells. Predominantly expressed in			
Antibody application notes  Positive control	For coating for ELISA, order Ab without BSA, Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Optimal dilution and staining procedure for a specific application should be determined by user, Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry  HL-60, MOLT-4 and MCF-7 cells. Predominantly expressed in ovary, prostate, testis and colon.			
Antibody application notes  Positive control  Shipping condition	For coating for ELISA, order Ab without BSA, Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Optimal dilution and staining procedure for a specific application should be determined by user, Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry  HL-60, MOLT-4 and MCF-7 cells. Predominantly expressed in ovary, prostate, testis and colon.  Room temperature  Store at 2 to 8 °C, Protect fluorescent conjugates from light,			
Antibody application notes  Positive control  Shipping condition Storage Conditions	For coating for ELISA, order Ab without BSA, Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Optimal dilution and staining procedure for a specific application should be determined by user, Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry  HL-60, MOLT-4 and MCF-7 cells. Predominantly expressed in ovary, prostate, testis and colon.  Room temperature  Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C			
Antibody application notes  Positive control  Shipping condition  Storage Conditions  Regulatory status  Antibody/conjugate	For coating for ELISA, order Ab without BSA, Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Optimal dilution and staining procedure for a specific application should be determined by user, Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry  HL-60, MOLT-4 and MCF-7 cells. Predominantly expressed in ovary, prostate, testis and colon.  Room temperature  Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C  For research use only (RUO)  Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified; 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, PIRFe: 1 mg/mL in PBS/0.05% BSA/0.05% azide, PIRFe: 1 mg/mL in PBS/0.05% BSA/0.05% azide, PIRFe: 1 mg/mL in			

Email: techsupport@biotium.com

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405	CF®405S Features
BNC88	CF®488A	490/515	488	GFP, FITC	CF®488A Features
BNC68	CF®568	562/583	532, 561	RFP, TRITC	CF®568 Features
BNC94	CF®594	593/614	561	Texas Red®	CF®594 Features
BNC40	CF®640R	642/662	633-640	Cy®5	CF®640R Features
BNC47	CF®647	650/665	633-640	Cy®5	CF®647 Features
BNC74	CF®740	742/767	633-685	775/50	CF®740 Features
BNCB	Biotin	N/A	N/A	N/A	
BNUB	Purified	N/A	N/A	N/A	
BNUM	Purified, BSA-free	N/A	N/A	N/A	

Alexa Fluor, Pacific Blue, Pacific Orange, and Texas Red are trademarks or registered trademarks of Thermo Fisher Scientific; Cy is a registered trademark of Cytiva; IRDye, LI-COR, and Odyssey are registered trademarks of LI-COR Bioscience.

This datasheet was generated on October 31, 2025 at 12:13:26 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/guanine-nucleotide-binding-protein-alpha-q-anaq-q-alpha-q-monoclonal-mouse-antibody-qnaq-2434/">https://biotium.com/product/guanine-nucleotide-binding-protein-alpha-q-anaq-q-alpha-q-monoclonal-mouse-antibody-qnaq-2434/</a>